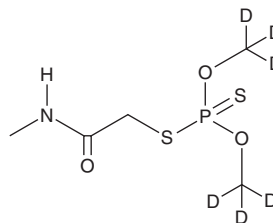


# PRODUCT INFORMATION



## Dimethoate-d<sub>6</sub> Item No. 39260

**CAS Registry No.:** 1219794-81-6  
**Formal Name:** phosphorodithioic acid, O,O-di(methyl-d<sub>3</sub>)  
S-[2-(methylamino)-2-oxoethyl] ester  
**MF:** C<sub>5</sub>H<sub>6</sub>D<sub>6</sub>NO<sub>3</sub>PS<sub>2</sub>  
**FW:** 235.3  
**Chemical Purity:** ≥98% (Dimethoate)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>6</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Dimethoate-d<sub>6</sub> is intended for use as an internal standard for the quantification of dimethoate (Item No. 24246) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Dimethoate-d<sub>6</sub> is supplied as a solid. A stock solution may be made by dissolving the dimethoate-d<sub>6</sub> in the solvent of choice, which should be purged with an inert gas. Dimethoate-d<sub>6</sub> is soluble in chloroform and DMSO.

### Description

Dimethoate is an organophosphate pesticide.<sup>1</sup> It is acaricidal against Kanzawa spider mites (LC<sub>50</sub> = 7 ppm in an aqueous suspension). Dimethoate (50-200 μM) reduces motility and viability and induces abnormal morphology of rat sperm.<sup>2</sup> It also increases production of malondialdehyde (MDA) and reduces superoxide dismutase (SOD), catalase (CAT), and glutathione peroxidase (GPX) activities in rat epididymal sperm. *In vivo*, dimethoate (28 mg/kg per day) decreases sperm count, motility, and viability and increases the percentage of morphologically abnormal sperm in rats.<sup>3</sup> Dimethoate also induces formation of carcinomas in the adrenal, thyroid, and pituitary glands of male and female rats as well as testicular atrophy, chronic renal disease, polyarteritis, and parathyroid hyperplasia in male rats.<sup>4</sup> Formulations containing dimethoate have been used as insecticides and miticides in aquatic food production and agriculture.

### References

1. Kuwahara, M. Insensitivity of the acetylcholinesterase from the organophosphate-resistant Kanzawa spider mite, *Tetranychus kanzawa* Kishida (Acarina: Tetranychidae) to organophosphorus and carbamate insecticides. *Appl. Ent. Zool.* **17(4)**, 486-493 (1982).
2. Ben Abdallah, F., Fetoui, H., Zribi, N., *et al.* Antioxidant supplementations *in vitro* improve rat sperm parameters and enhance antioxidant enzyme activities against dimethoate-induced sperm damages. *Andrologia* **44(Suppl 1)**, 272-279 (2012).
3. Abdallah, F.B., Slima, A.B., Dammak, I., *et al.* Comparative effects of dimethoate and deltamethrin on reproductive system in male mice. *Andrologia* **42(3)**, 182-186 (2010).
4. Reuber, M.D. Carcinogenicity of dimethoate. *Environ. Res.* **34(2)**, 193-211 (1984).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/10/2023

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM